

# PRODUCT INFORMATION



## SARS-CoV-2 Spike Glycoprotein S1 Subunit

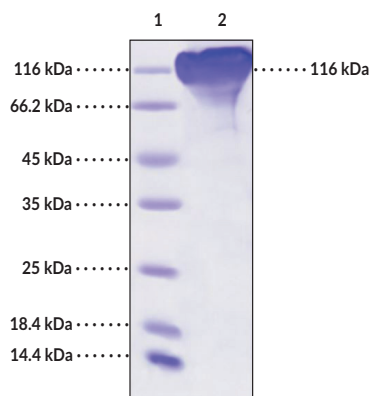
Item No. 31814

### Overview and Properties

<b>Synonyms:</b>	COVID-19 Surface Glycoprotein S1 Subunit, 2019-nCoV Surface Glycoprotein S1 Subunit, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein S1 Subunit
<b>Source:</b>	Active recombinant C-terminal mouse IgG1 Fc-tagged SARS-CoV-2 spike glycoprotein S1 subunit expressed in HEK293 cells
<b>Amino Acids:</b>	16-685
<b>Uniprot No.:</b>	PODTC2
<b>Molecular Weight:</b>	101.4 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥90% estimated by SDS-PAGE
<b>Supplied in:</b>	Lyophilized from sterile PBS, pH 7.4
<b>Endotoxin Testing:</b>	<1.0 EU/g, determined by the LAL endotoxin assay
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml
<b>Bioactivity:</b>	See figures for details

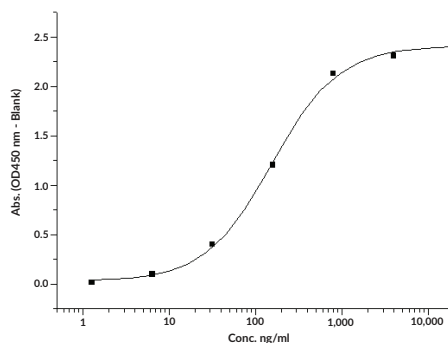
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: SARS-CoV-2 Spike Glycoprotein S1 Subunit

**SDS-PAGE Analysis of SARS-CoV-2 Spike Glycoprotein S1 Subunit.**  
This protein has a calculated molecular weight of 101.4 kDa. It has an apparent molecular weight of approximately 116 kDa by SDS-PAGE under reducing conditions due to glycosylation.



**SARS-CoV-2 Spike Glycoprotein S1 Subunit Activity in a Functional ELISA.** Immobilized human ACE2 protein (His-tag) at 2 µg/ml (100 µl/well) can bind SARS-CoV-2 Surface Glycoprotein S1 Subunit. The EC<sub>50</sub> value for this is typically 90-240 ng/ml.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/09/2021

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) surface glycoprotein, also known as the spike glycoprotein, is encoded by the *S* gene in SARS-CoV-2 RNA.<sup>1</sup> SARS-CoV-2 is a member of the *Betacoronavirus* genus of viruses that has an approximately 79% sequence identity with SARS-CoV.<sup>2,3</sup> The spike protein of SARS-CoV-2 and the related viruses SARS-CoV and Middle East respiratory syndrome coronavirus (MERS-CoV) is a transmembrane glycoprotein that assembles into homotrimers on the virus surface and is comprised of an N-terminal S1 subunit, which contains the receptor binding domain (RBD), and a C-terminal S2 subunit, which facilitates fusion between viral and host cell membranes.<sup>4-6</sup> The 193-amino acid RBD of the SARS-CoV spike protein is a target for neutralizing antibodies.<sup>5,7</sup> The SARS-CoV-2 RBD, which spans amino acid residues 329 to 521, is 73% identical to that of SARS-CoV and can bind to human angiotensin-converting enzyme 2 (ACE2), which is the host cell surface receptor for both SARS-CoV and SARS-CoV-2.<sup>4-7</sup> SARS-CoV-2 is the causative agent of COVID-19, a primarily respiratory illness characterized by fever, cough, and shortness of breath that can lead to life-threatening complications.<sup>6,8,9</sup> Cayman's SARS-CoV-2 Surface Glycoprotein S1 Subunit can be used for ELISA. This protein is a disulfide-linked homodimer. The reduced monomer, comprised of the SARS-CoV-2 surface glycoprotein S1 subunit (amino acids 16-685) fused to mouse IgG1 Fc at its C-terminus, consists of 904 amino acids and has a calculated molecular weight of 101.4 kDa. As a result of glycosylation, the monomer migrates at approximately 116 kDa by SDS-PAGE under reducing conditions.

## References

---

1. Kandeel, M., Ibrahim, A., Fayez, M., *et al.* From SARS and MERS CoVs to SARS-CoV-2: Moving toward more biased codon usage in viral structural and nonstructural genes. *J. Med. Virol.* **92(6)**, 660-666 (2020).
2. Lu, R., Zhao, X., Li, J., *et al.* Genomic characterisation and epidemiology of 2019 novel coronavirus: Implications for virus origins and receptor binding. *Lancet* **395(10224)**, 565-574 (2020).
3. Meo, S.A., Alhowikan, A.M., Al-Khlaiwi, T., *et al.* Novel coronavirus 2019-nCoV: Prevalence, biological and clinical characteristics comparison with SARS-CoV and MERS-CoV. *Eur. Rev. Med. Pharmacol. Sci.* **24(4)**, 2012-2019 (2020).
4. Liu, Z., Xiao, X., Wei, X., *et al.* Composition and divergence of coronavirus spike proteins and host ACE2 receptors predict potential intermediate hosts of SARS-CoV-2. *J. Med. Virol.* **92(6)**, 595-601 (2020).
5. He, Y., Zhou, Y., Liu, S., *et al.* Receptor-binding domain of SARS-CoV spike protein induces highly potent neutralizing antibodies: Implication for developing subunit vaccine. *Biochem. Biophys. Res. Commun.* **324(2)**, 773-781 (2004).
6. Walls, A.C., Park, Y.-J., Tortorici, M.A., *et al.* Structure, function, and antigenicity of the SARS-CoV-2 spike glycoprotein. *Cell* **181(2)**, 281-292 (2020).
7. Tian, X., Li, C., Huang, A., *et al.* Potent binding of 2019 novel coronavirus spike protein by a SARS coronavirus-specific human monoclonal antibody. *Emerg. Microbes Infect.* **9(1)**, 382-385 (2020).
8. Klok, F.A., Kruip, M.J.H.A., van der Meer, N.J.M., *et al.* Incidence of thrombotic complications in critically ill ICU patients with COVID-19. *Thromb. Res.* **191**, 145-147 (2020).
9. Yang, F., Shi, S., Zhu, J., *et al.* Analysis of 92 deceased patients with COVID-19. *J. Med. Virol.* (2020).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM